

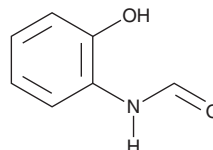
# PRODUCT INFORMATION



## N-formyl-2-Aminophenol

Item No. 36802

**CAS Registry No.:** 2843-27-8  
**Formal Name:** N-(2-hydroxyphenyl)-formamide  
**Synonyms:** 2'-Hydroxyformanilide, NSC 93900  
**MF:** C<sub>7</sub>H<sub>7</sub>NO<sub>2</sub>  
**FW:** 137.1  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 245 nm  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

N-formyl-2-Aminophenol is supplied as a solid. A stock solution may be made by dissolving the N-formyl-2-aminophenol in the solvent of choice, which should be purged with an inert gas. N-formyl-2-Aminophenol is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of N-formyl-2-aminophenol in these solvents is approximately 1 mg/ml.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of N-formyl-2-aminophenol can be prepared by directly dissolving the solid in aqueous buffers. The solubility of N-formyl-2-aminophenol in PBS (pH 7.2) is approximately 0.16 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

N-formyl-2-Aminophenol is a bacterial secondary metabolite that has been found in *P. chrysogenum* and has antioxidant activity.<sup>1,2</sup> It scavenges DPPH (Item No. 14805) radicals in a cell-free assay with an IC<sub>50</sub> value of 3.23 μg/ml.<sup>2</sup> N-formyl-2-Aminophenol has also been used in the synthesis of compounds with antibacterial and antifungal activities.<sup>3</sup>

### References

1. Frisvad, J.C., Smedsgaard, J., Larsen, T.O., *et al.* Mycotoxins, drugs and other extrolites produced by species in *Penicillium* subgenus *Penicillium*. *Stud. Mycol.* **49**, 201-241 (2004).
2. Deepali, A., Akansha, A., Anamika, B., *et al.* Hydroxylamine hydrochloride as an effective catalyst for formamide derivative synthesis and their DPPH scavenging activity. *Res. J. Chem. Sci.* **4(10)**, 54-57 (2014).
3. Gupta, S., Verma, P., and Singh, V. Synthesis and antimicrobial study of 2-amino-imidazole derivatives. *Indian J. Chem.* **57B**, 679-686 (2018).

#### WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

#### SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 12/21/2022

#### CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA

**PHONE:** [800] 364-9897  
[734] 971-3335

**FAX:** [734] 971-3640

CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM